

File :Y:\Forensics\11011-0447669-2.D
Operator : PDK
Acquired : 26 Jun 2013 13:10 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: Pentedrone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 1

